

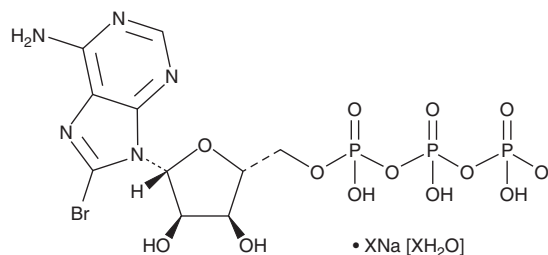
# PRODUCT INFORMATION



## 8-Bromoadenosine 5'-triphosphate (sodium salt hydrate)

Item No. 30573

<b>Formal Name:</b>	8-bromo-adenosine 5'-(tetrahydrogen triphosphate), sodium salt hydrate
<b>Synonym:</b>	8-bromo ATP
<b>MF:</b>	$C_{10}H_{15}BrN_5O_{13}P_3 \cdot XNa [XH_2O]$
<b>FW:</b>	586.1
<b>Purity:</b>	≥95%
<b>UV/Vis.:</b>	$\lambda_{max}$ : 264 nm
<b>Supplied as:</b>	A solid
<b>Storage:</b>	-20°C
<b>Stability:</b>	≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

8-Bromoadenosine 5'-triphosphate (8-Br-ATP) (sodium salt hydrate) is supplied as a solid. A stock solution may be made by dissolving the 8-Br-ATP (sodium salt hydrate) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

8-Br-ATP is an analog of ATP (Item No. 14498) and an agonist of the purinergic P2X receptor.<sup>1</sup> It induces contraction of isolated guinea pig bladder strips with a potency that is 0.19-fold that of ATP. 8-Br-ATP has been used in the determination of the nucleotide substrate specificity of yeast poly(A) polymerase.<sup>2</sup> It is cytotoxic to multiple myeloma cells ( $IC_{50} = 23.1 \mu M$ ).<sup>3</sup>

### References

1. Howson, W., Taylor, E.M., Parsons, M.E., *et al.* Synthesis and biological evaluation of ATP analogues acting at putative purinergic P<sub>2X</sub>-receptors (on the guinea pig bladder). *Eur. J. Med. Chem.* **23(5)**, 433-439 (1988).
2. Chen, L.S. and Sheppard, T.L. Chain termination and inhibition of *Saccharomyces cerevisiae* poly(A) polymerase by C-8-modified ATP analogs. *J. Biol. Chem.* **279(39)**, 40405-40411 (2004).
3. Wang, L. and MacDonald, R.C. Cationic phospholiposomes: Efficient delivery vehicles of anticancer derivatives of ATP to multiple myeloma cells. *J. Liposome Res.* **21(4)**, 306-314 (2011).

#### WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

#### SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/12/2022

#### CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897  
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM